



In Vitro Skin Corrosion Test: Reproducibility Over Time and Optimized Methodology for Testing Chemicals Interfering with the MTT endpoint

Helena Kandarova¹, Hans Raabe², Gregory Chua¹, Mitch Klausner¹, Joseph Kubilus¹, Patrick Hayden¹, Seyoum Ayeahunie¹, Yulia Kaluzhny¹, Rodger Curren² and John Sheasgreen¹

¹MatTek Corporation, Ashland, MA

²Institute for In Vitro Science, Inc. Gaithersburg, MD

The potential for chemicals to cause skin effects such as corrosion is a concern of industrial toxicologists in their assessments of possible worker and consumer safety issues. Moreover, the U.S. Department of Transportation (DOT), the new regulation on the Registration, Evaluation and Authorization of Chemicals (REACH), and other international and national regulatory agencies require that substances should be labeled as to skin corrosivity potential.

In the past, skin corrosion assessments were based on tests involving topical application of test substances to the skin of rabbits. However, based on two ECVAM Validation studies performed during 1996-2000 (Fentem et al., *Toxicology in Vitro* 12, 483-524, 1998; Liebsch et al., *ATLA* 28, 371-401, 2000) with two reconstructed human skin models (EpiDerm and EPIKIN), the OECD approved use of skin models as regulatory accepted methods (OECD TG 431) replacing the in vivo test.

In the present study, the EpiDerm skin corrosion test was repeated with the commercially available test substances previously used in both ECVAM validation studies. The aim was to demonstrate the long term reproducibility and reliability of the EpiDerm model and the method, as required by regulators (Rispin et al., *Regul. Toxicol. Pharmacol.* 45 (2), 97-103, 2006). The data obtained show very good correlation over a period of 7 years. We also conducted a formalized evaluation of a procedural modification proposed several years ago to correctly predict the corrosive potential of materials that interfere with the MTT endpoint normally used in the method. An attempt was made to optimize this method and demonstrate with several chemicals (sometimes reported to be false negatives) the predictive capacity of the EpiDerm corrosivity assay when this procedural modification is included. The presentation will summarize data from the above mentioned studies and describe optimized methods for assessing materials which interfere with the MTT endpoint.

Keywords: EpiDerm, skin corrosion, REACH, MTT endpoint, regulatory toxicology

Session: EU-chemicals policy (REACH)

To be presented at the Linz 2007 – 14th Congress on Alternatives to Animal Testing, September 28-30, 2007 at the University of Linz, Austria.